

# PRODUCT INFORMATION



**SJ000291942**

Item No. 29904

**CAS Registry No.:** 425613-09-8  
**Formal Name:** 2-(4-ethylphenoxy)-N-(4-fluoro-3-nitrophenyl)-acetamide

**MF:** C<sub>16</sub>H<sub>15</sub>FN<sub>2</sub>O<sub>4</sub>

**FW:** 318.3

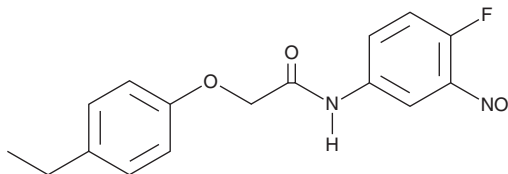
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 235 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

SJ000291942 is supplied as a crystalline solid. A stock solution may be made by dissolving the SJ000291942 in the solvent of choice, which should be purged with an inert gas. SJ000291942 is soluble in DMSO and dimethyl formamide. The solubility of SJ000291942 in these solvents is 30 mg/ml, respectively.

## Description

SJ000291942 is a bone morphogenic protein (BMP) signaling activator.<sup>1</sup> It activates BMP4 in C33A-2D2 human cervical carcinoma clonal reporter cells (EC<sub>50</sub> = ≤1 μM). SJ000291942 (0.1-50 μM) induces ventralization, as well as increases expression of the BMP target genes *bmp2b* and *szl* and the cooperatively-induced BMP target genes *vent* and *vox* in zebrafish embryos. It also induces BMP-dependent osteoblastic differentiation of C2C12 murine myoblasts.

## Reference

1. Genthe, J.R., Min, J., Farmer, D.M., *et al.* Ventromorphins: A new class of small molecule activators of the canonical BMP signaling pathway. *ACS Chem. Biol.* **12(9)**, 2436-2447 (2017).

### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/14/2022

## CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM